

# Splitting Solution

## ***Animal Product-Free***

Catalog Number: R-008-25  
25 ml

Catalog Number: R-008-3  
3 ml

## **Instructions for storage and use**

### **Product Description**

Splitting Solution is an animal product-free dispase solution (1X, 25 units/ml) for separating epidermis from dermis.

### **Intended Use**

Splitting Solution is intended for use as one component for the routine isolation from tissue of normal human epidermal and dermal cells in an animal product-free environment. Splitting Solution in the 3 ml size (cat. # R-008-3) is one component in Cascade Biologics' Keratinocyte Primary Isolation Kit (cat. # R-014-K).

***This product is for research use only. Not for use in animals, humans, or diagnostic procedures.***

***Caution: If handled improperly, some components of this product may present a health hazard. Take appropriate precautions when handling this product, including the wearing of protective clothing and eyewear. Dispose of properly.***

### **Storage and Stability**

Splitting Solution is stored at -20° C at our facility and is shipped on dry ice. Upon receipt, Splitting Solution should be stored at -20° C in a freezer that is not self-defrosting. When stored at -20° C, the product is stable until the expiration date on the label.

After long-term storage at -20° C, Splitting Solution may contain a small amount of precipitate. This precipitate will not affect the performance of the product.

### **Instructions for Use**

*For detailed instructions on isolating normal human keratinocytes from tissue using Splitting Solution, please refer to the product instructions that accompany our Keratinocyte Primary Isolation Kit (cat. # R-014-K), which can be found at: [www.cascadebio.com](http://www.cascadebio.com).*

## **Terms and Conditions of Sale**

Cascade Biologics, Inc. (hereinafter, CBI) warrants that its products will perform according to the information provided in various publications that it distributes, and as described herein for the intended shelf life of the product when stored under the conditions prescribed by CBI. If any product does not perform according to the published information provided by us, CBI will replace the product free of charge to the original customer. The remedy of product replacement shall be the customers' sole and exclusive remedy for defective product, unless CBI is unable to deliver replacement product, in which case CBI shall reimburse the customer for the purchase price of the defective product. Customer understands that the foregoing limited warranty is in lieu of all other warranties and CBI hereby disclaims all other warranties including, but not limited to, implied warranty of merchantability or fitness for adequacy for any particular purpose or use. CBI shall not be liable to customer or to any party claiming through customer for any incidental or consequential damages, including, but not limited to any lost profits, lost savings, or lost business, whether arising out of contract, tort, or otherwise. Customer acknowledges that products are intended for research use only and are not to be used in animals, humans, or diagnostic procedures.